

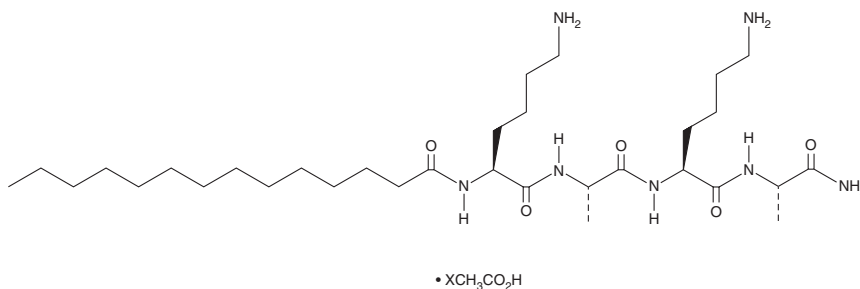
PRODUCT INFORMATION



Myristoyl Tetrapeptide-12 (acetate)

Item No. 45604

Formal Name: N²-(1-oxotetradecyl)-L-lysyl-L-alanyl-L-lysyl-L-alaninamide, acetate
Synonym: Myr-Lys-Ala-Lys-Ala-NH₂
Peptide Sequence: Myr-KAKA-NH₂
MF: C₃₂H₆₃N₇O₅ • XC₂H₄O₂
FW: 625.9
Purity: ≥98%
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

Myristoyl tetrapeptide-12 (acetate) is supplied as a solid. Aqueous solutions of myristoyl tetrapeptide-12 (acetate) can be prepared by directly dissolving the solid in aqueous buffers. Myristoyl tetrapeptide-12 (acetate) is sparingly soluble (1-10 mg/ml) in PBS (pH 7.2). We do not recommend storing the aqueous solution for more than one day.

Description

Myristoyl tetrapeptide-12 is a polymeric acylated biosurfactant.¹ It has a critical micelle concentration of less than 200 ppm in aqueous minimal essential media (MEM) and is active against *E. coli*, *C. albicans*, *P. aeruginosa*, and *S. aureus* (MICs = ≤100 ppm). Formulations containing myristoyl tetrapeptide-12 have been used in cosmetic products.

Reference

1. Owen, D., and Fan, L. Polymeric biosurfactants. *Therapeutic Peptides, Inc.* **WO 2007/143006 A2** (2007).

WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the [complete](#) Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 05/12/2026

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897

[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM